

Research Associate I

Luceome Biotechnologies, an emerging biotechnology company, provides unique assay solutions to clients for selectivity profiling of drugs against specific targets. Luceome's assays are based on a technology platform that utilizes luciferase fragment complementation, which allows identification of target and off-target effects of drug candidates with high sensitivity and minimal background interference. For more information, please visit www.luceome.com.

Position Overview

Luceome is seeking research oriented, motivated person to contribute to its *KinaseSeeker*[™] program for profiling drug candidates against kinases. The internship candidate will assist in building a large panel of kinases, validating the kinases in split-luciferase based assays, and evaluating test compounds in these assays. Candidates should preferably have research experience in molecular biology techniques, especially cloning, PCR and gel electrophoresis; and biochemical assays. Familiarity with kinases is not required but will be regarded as a plus. Preference will be given to candidates with high analytical ability and excellent communication skills.

Qualifications:

- B.S. or M.S in Biochemistry or Molecular Biology. Candidate should have worked in a laboratory setting for a minimum of 2 years
- Strong track record of scientific achievement
- Hands-on experience in cloning, PCR, gel electrophoresis and biochemical assays
- Willingness to work in a dynamic and changing work environment
- Excellent written and verbal skills

Interested candidates are encouraged to send their CV with cover letter to info@luceome.com.